Maxwell® RSC Tissue DNA Kit

Instructions for Use of Product AS1610.



Preparing Tissue Samples for DNA Purification

This Quick Protocol provides instructions for use of the Maxwell® RSC Tissue DNA Kit with Maxwell® Instruments to purify genomic DNA from up to 50mg of tissue. For detailed instructions, including information on instrument setup and troubleshooting, please refer to the *Maxwell® RSC Tissue DNA Kit Technical Manual* #TM476, available at: **www.promega.com/protocols/**

Materials to be Supplied by the User

- pipettors and pipette tips for sample transfer
- Potter-Elvehjem PTFE pestle and glass tube (recommended) or Disposable Pellet Pestles with Tube (Fisher Cat.# 12-141-386) The total yield of genomic DNA from tissue samples is variable and depends on the type of tissue being used.

Note: Some tissues will require vigorous disruption techniques. For these tissues, a Potter-Elvehjem homogenizer is strongly recommended. Use of tissue samples that are not fully disrupted may result in poor purification performance and tissue may transfer into the elution tube.

- 1. Weigh out up to 50mg of tissue sample and place in tube.
- 2. Add 80µl of TE buffer to the sample in the tube.
- 3. Using the pestle, thoroughly disrupt and homogenize the tissue sample.

Maxwell® Automated DNA Purification

Cartridge Preparation

- 1. Place the cartridge to be used in the deck tray with well #1 (the largest well in the cartridge) facing away from the elution position, which is the numbered side of the tray.
- 2. Press down on the cartridge to snap it into position. Carefully peel back the seal so that all plastic comes off the top of the cartridge. Ensure that all sealing tape and any residual adhesive are removed before placing the cartridge in the instrument.
- 3. Transfer homogenized tissue sample to well #1 (the largest well) of each cartridge and thoroughly mix the tissue samples into the lysis buffer by pipetting at least 10 times.
- 4. Place one plunger in well #8 of each cartridge. Well #8 is the well closest to the elution tube.
- 5. Place an empty elution tube into the elution tube position for each cartridge. Add 100µl of Elution Buffer to the bottom of each elution tube.

Note: Use only the Elution Tubes (0.5ml) provided with the kit; other tubes may be incompatible with supported Maxwell[®] Instruments.

Instrument Run on the Maxwell® Instruments

Follow the instrument run instructions in the Maxwell® RSC Tissue DNA Kit Technical Manual #TM476.

Maxwell is a registered trademark of Promega Corporation.

Additional protocol information is in Technical Manual #TM476, available online at: www.promega.com

